CD4 Monoclonal Mouse Antibody (CD4/1604)



Product Description

The immunogen for this Mab is derived from the extracellular domain of CD4, a membrane glycoprotein of T lymphocytes that interacts with major histocompatibility complex class II antigens and is also a receptor for the human immunodeficiency virus. This protein is expressed not only in T lymphocytes, but also in B cells, macrophages, and granulocytes. It is also expressed in specific regions of the brain. The protein functions to initiate or augment the early phase of T-cell activation, and may function as an important mediator of indirect neuronal damage in infectious and immune-mediated diseases of the central nervous system. The majority of peripheral T-cell lymphomas are derived from the T-helper/regulatory cell subset so that most mature T-cell neoplasms are CD4 /CD8-. Anti-CD4 is used in the immunohistochemical staining of lymphoproliferative disorders to evaluate tumors with CD4 aberrant expression.

Also view our <u>CD4 Monoclonal Mouse Biotium Choice Antibody (SK3)</u>. The antibody belongs to the Biotium Choice list of select antibodies that have been developed and validated in-house for optimal performance. The antibody has been validated in flow cytometry.

Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the CF® Dye Brochure for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors.

Stock status: Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email order@biotium.com to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 1604, CD4 Monoclonal Mouse Antibody (CD4/1604)

Droduct attributor

Cell/tissue expression

Tumor expression

Call us: 800-304-5357

| Product attributes | | | | |
|---------------------------------------|--|--|--|--|
| Antibody number | #1604 | | | |
| Antibody reactivity (target) | CD4 | | | |
| Antibody type | Primary | | | |
| Host species | Mouse | | | |
| Clonality | Monoclonal | | | |
| Clone | CD4/1604 | | | |
| Isotype | IgG2b, kappa | | | |
| Molecular weight | 55 kDa | | | |
| Synonyms | L3T4; Leu3; Ly-4; Lymphocyte antigen CD4; p55; T cell antiger T4/LEU3; T cell differentiation antigen L3T4; T-cell surface antigen T4/Leu-3; T-cell surface glycoprotein CD4 | | | |
| Human gene symbol | CD4 | | | |
| Entrez gene ID | 920 | | | |
| SwissProt | P01730 | | | |
| Unigene | 631659 | | | |
| Immunogen | Recombinant human CD4 protein extracellular domain fragment (around aa 245-392) (exact sequence is proprietary) | | | |
| Antibody target cellular localization | Plasma membrane | | | |
| Verified antibody applications | Flow (surface) (verified) | | | |
| Species reactivity | Human | | | |
| Antibody application notes | ELISA: For coating, order Ab without BSA; Optimal dilution for a specific application should be determined., Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody | | | |
| Positive control | Human peripheral blood mononuclear cells (PBMCs), CCRF-CEM or HL-60. Lymph node or tonsil. | | | |
| Shipping condition | Room temperature | | | |
| Storage Conditions | Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C | | | |
| Regulatory status | For research use only (RUO) | | | |
| Antibody/conjugate formulation | Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/ml in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL PBS without azide | | | |
| Shelf life | Guaranteed for at least 24 months from date of receipt when stored as recommended | | | |
| Product origin | roduct origin Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot. | | | |
| | | | | |

Lymphocytes, Myeloid cells

Leukemia/lymphoma

| Antibody # prefix | Conjugation | Ex/Em (nm) | Laser line | Detection channel | Dye Features |
|-------------------|-----------------------|------------|------------|-----------------------------|------------------|
| BNC04 | CF®405S | 404/431 | 405 | DAPI (microscopy), AF405 | CF®405S Features |
| BNC88 | CF®488A | 490/515 | 488 | GFP, FITC | CF®488A Features |
| BNC68 | CF®568 | 562/583 | 532, 561 | RFP, TRITC | CF®568 Features |
| BNC94 | CF®594 | 593/614 | 561 | Texas Red® | CF®594 Features |
| BNC40 | CF®640R | 642/662 | 633-640 | Cy®5 | CF®640R Features |
| BNC47 | CF®647 | 650/665 | 633-640 | Cy®5 | CF®647 Features |
| BNC74 | CF®740 | 742/767 | 633-685 | 775/50 | CF®740 Features |
| BNCB | Biotin | N/A | N/A | N/A | |
| BNUB | Purified | N/A | N/A | N/A | |
| BNUM | Purified, BSA-free | N/A | N/A | N/A | |

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, and Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on November 30, 2025 at 04:25:10 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/cd4-monoclonal-mouse-antibody-cd4-1604/